PRODUCT INFORMATION



Phenazepam-d₄ Item No. 16198

CAS Registry No.: 1184980-42-4

Formal Name: 7-bromo-5-(6-chlorophenyl-

2,3,4,5-d₄)-1,3-dihydro-2H-1,4-

benzodiazepin-2-one

Synonym: Fenazepam-d₄ MF: C₁₅H₆BrClD₄N₂O

FW: 353.6 **Purity:** ≥98% Supplied as: A neat solid Storage: -20°C Stability: ≥4 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Phenazepam-d₄ (Item No. 16198) is an analytical reference material intended for use as an internal standard for the quantification of phenazepam (Item Nos. 9000849 | ISO60172) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Phenazepam is categorized as a benzodiazepine. It has sedative and hypnotic properties and has been detected in illicit tablets.² This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

- 1. Kalinina, T.S., Garibova, T.L., and Voronina, T.A. Descriminative effects of phenazepam and gidazepam in rats: Comparison with other GABA-related drugs. Pharmacol. Biochem. Behav. 64(2), 397-401 (1999).
- Oyemade, A. New uncontrolled benzodiazepine, phenazepam, emerging drug of abuse. Innov. Clin. Neurosci. 9(9), 10-13 (2012).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the material can be found on our website.

Copyright Cayman Chemical Company, 07/10/2019

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD ANN ARBOR, MI 48108 · USA PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640 CUSTSERV@CAYMANCHEM.COM WWW.**CAYMANCHEM**.COM